

Recombinant Rat Interleukin-6 (IL-6)








Publication Number MAN0004413





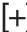

Rev. 1.00

Catalog Number:	PRC0064	PRC0065	PRC0061	PRC0063
Quantity:	10 µg	25 µg	100 µg	1 mg
Lot Number:	See product label.			
Molecular Weight:	22 kDa			
Purity:	>95% as determined by SDS-PAGE analysis.			
Biological Activity:	ED ₅₀ = 0.03–0.1 ng/mL. The biological activity is determined by measuring the dose dependent proliferation of IL-6 dependent B9 cells. A concentration range of 0.1–10.0 ng/mL is effective for most <i>in vitro</i> applications. The optimal concentration for each specific application should be determined by an initial dose response assay.			
Formulation:	Lyophilized, carrier free.			
Sterility:	Filtered prior to lyophilization through a 0.22 micron sterile filter.			
Endotoxin:	<0.1 ng/µg			
Production:	Recombinant rat IL-6 is produced in <i>E. coli</i> and purified via sequential chromatography.			
Reconstitution Recommendation:	We recommend that this vial be briefly centrifuged prior to opening to bring the contents to the bottom. Lyophilized rIL-6 should be reconstituted in 100 mM acetic acid to 1.0 mg/mL and incubated for 30 minutes at room temperature to regain full activity. Stock solutions should be apportioned into working aliquots and stored at ≤ –20°C. Further dilutions should be made in low endotoxin medium or buffered solution with FCS or tissue culture grade BSA.			
Suggested Working Dilutions:	The optimal concentration should be determined for each specific application.			
Storage:	Lyophilized rIL-6 should be stored at 2°C to 8°C, preferably desiccated. Store reconstituted rIL-6 at ≤ –20°C (not in a frost-free freezer). Keep freeze–thaw cycles to a minimum.			
Expiration Date:	Expires one year from date of receipt when stored as instructed.			
References:	<p>Northemann, W., T.A. Braciak, M. Hattori, F. Lee, and G.H. Fey (1989) Structure of the rat interleukin 6 gene and its expression in macrophage-derived cells. <i>J. Biol. Chem.</i> 264:16072–16082.</p> <p>Aarden, L.A., E.R. De Groot, O.L. Schaap, and P.M. Landsorp (1987) Production of hybridoma growth factor by human monocytes. <i>Eur. J. Immunol.</i> 17:1411–1416.</p> <p>Hirano, T. (1994) In: <i>The Cytokine Handbook</i>, 2nd Edition (ed. A. Thompson) Academic Press Limited, San Diego, CA, pp. 145–168.</p> <p>Rees, G.S., C. Ball, H.L. Ward, C.K. Gee, G. Tarrant, Y. Mistry, S. Poole, and A.F. Bristow (1999) Rat interleukin 6: expression in recombinant <i>Escherichia coli</i>, purification and development of a novel ELISA. <i>Cytokine</i> 11:95–103.</p> <p>Ikeda, Y., A. Sugawara, Y. Taniyama, A. Uruno, K. Igarashi, S. Arima, S. Ito, and K. Takeuchi (2000) Suppression of rat thromboxane synthase gene transcription by peroxisome proliferator-activated receptor γ in macrophages via an interaction with NRF2. <i>J. Biol. Chem.</i> 275(42):33142–33150.</p>			

Explanation of Symbols

The symbols present on the product label are explained below:

Symbol	Description
	Catalog Number
	Research Use Only
	Use by
	Manufacturer
	Without, does not contain
	Protect from light
	Directs the user to consult instructions for use (IFU), accompanying the product.

Symbol	Description
	Batch code
	In vitro diagnostic medical device
	Temperature limitation
	European Community authorized representative
	With, contains
	Consult accompanying documents

Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

For Research Use Only. Caution: Not for human or animal therapeutic or diagnostic use.

Manufactured under ISO 13485 Quality Standard

Manufacturing site: 7335 Executive Way | Frederick, MD 21704 | Toll Free in USA 800.955.6288

© 2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

For support visit www.lifetechnologies.com/support or email techsupport@lifetech.com

www.lifetechnologies.com

Revision Date 24 May 2011

